

Certificate of Analysis

Human IL-6, human cell-expressed (HCE)

Catalog No: 11605-2

Lot No: H00052

Expiration: One year from date of receipt

Size: 10 µg

Description: HumanKine® Recombinant Human IL-6

EC50: Typically 0.2 to 1 ng/ml

Formulation: Lyophilized from a 0.2 µm filtered solution of PBS

Endotoxin: < 1 EU/µg

Molecular Weight: 26 and 30 kDa, monomer, glycosylated

Purity: > 95%

Purity Confirmation: SDS-PAGE stained with Coomassie blue

Source: Human Cell Expressed (HCE) in HEK-293 cells in the absence of animal-derived products

Reconstitution: Briefly centrifuge the vial before opening. It is recommended to reconstitute the protein in sterile PBS containing 0.1% recombinant human serum albumin, low endotoxin (please see product catalog #11600 HSA).

Assay Used to Measure Bioactivity: IL-6 specific activity was determined by the dose-dependent proliferation of mouse 7TD1 cells.

References:

1. Kitamura, T., *et al.* (1989) *J. Cell Physiol.* 140, 323-334.
2. Naugler, W.E., *et al.* (2008) *Trends Mol Med* doi:10.1016/j.molmed.2007.12.007

Product Information:

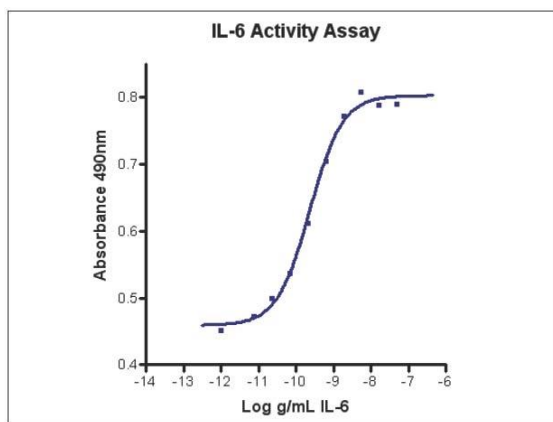


Figure 1. Human IL-6 Activity Assay

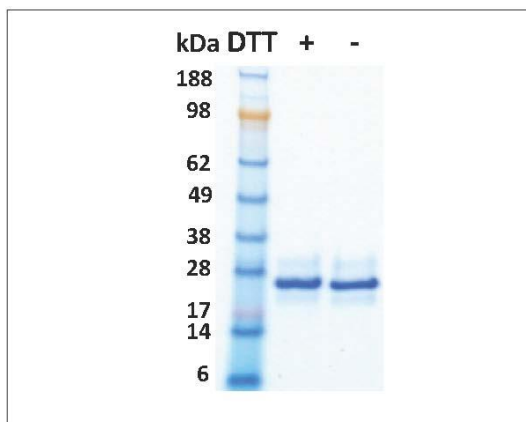


Figure 2. Purity Confirmation on SDS-PAGE



Shipping Conditions: Room Temperature, 22-25°C

Physical State of Product During Shipping: Lyophilized

Storage Conditions/Comments: Lyophilized products are stable for one year at -20°C to -80°C from date of receipt. Upon reconstitution, the cytokine can be stored at 4°C for short term only (< 1 week), or at -20°C to -80°C in aliquots for long term. Avoid repeated freeze-thaw cycles. Refreezing should be performed on dry ice or in a dry ice/alcohol bath. For more information on protein handling, visit our Resource Library at www.pblsaysci.com

Authorization

Released by: _____

A handwritten signature in black ink, appearing to be 'L. J. ...', written over a horizontal line.

Date: March 15, 2016

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL Assay Science. For research use only. Not for diagnostic or clinical use in, or administration to, humans. Not for resale in original or any modified form, including inclusion in a kit, for any purpose. Not for use in the preparation of any commercial product.